EUROPE

LIEGE SCIENCE PARK • 4102 Seraing • BELGIUM • Tel.: +32 4 372 74 00 • Fax: +32 4 372 75 00 Toll-free: +800 666 00 123 • info@eurogentec.com • www.eurogentec.com

NORTH AMERICA

ANASPEC - 34801 Campus Drive • Fremont, CA 94555 • USA • Tel.: +1-510-791-9560 Toll-free: +1 800-452-5530 • Fax: +1 510-791-9572 • service@anaspec.com



Technical Data Sheet

Mu-MLV Reverse Transcriptase

ME-0125-400

Eurogentec products are sold for research or laboratory use only and are not to be administrated to humans or used for medical diagnostics.

Applications

Mu-MLV reverse transcriptase is able to synthesize a DNA strand complementary to an RNA template in the presence of a primer. This enzyme is particularly suited for long fragments.

Batch details

Units per vial: ME-0125-400 40000 units

Concentration: 200 U/µl

Source

The enzyme is isolated from a recombinant E.coli clone over expressing the enzyme.

Description

Mu-MLV reverse transcriptase has lower RNAse H activity than AMV reverse transcriptase, which is an advantage when synthesizing cDNAs from long mRNAs. Mu-MLV lacks the $3' \rightarrow 5'$ exonuclease activity. It is recommended to use 10 U/µg RNA for optimum reverse transcription.

Although the enzyme shows low activity at 42 $^{\circ}$ C, it is quite stable at 37 $^{\circ}$ C.

Package contents

Reagent	Description	
Mu-MLV reverse transcriptase	ME-0125-400	
5x cDNA buffer (clear cap)	250 mM Tris-HCl (pH 8.3),	
	375 mM KCl, 50 mM DTT,	
	15 mM MgCl ₂	

Shipping conditions

Shipping at 4 °C.

Storage conditions

Storage at -20 °C is recommended.

Storage buffer

50 mM Tris-HCl pH 8.3 (4 °C), 1 mM EDTA, 0.1 % Triton X-100, 0.1 M NaCl, 5 mM DTT, 50 % (ν V) Glycerol.

Analysis conditions

50 mM Tris-HCl (pH 8.3), 75 mM KCl, 3 mM MgCl $_{\rm 2}$, 0.5 mM dTTP, 10 mM DTT, 2.5 μ M polyA-oligodT 12-18, incubate at 37 °C.

Unit definition

One unit is the amount of enzyme that incorporates 1 nmole of dTTP into acid - insoluble form in 10 minutes at 37 °C using polyA-oligodT 12-18 as substrate.

Reaction procedure

- Thaw all required reagents except for the Mu-MLV reverse transcriptase, which should be kept in the freezer until use. Mix all reagents by inversion and spin them down prior to pipeting. Note: To correct for dispensing losses prepare an excess of reaction mix. A negative control containing no RNA template should always be included.
- 2. Prepare the RT reaction mix by adding the following components to a nuclease-free 0.2 ml thermocycler tube:
 - x μl oligo d(T)12-18, or random primers, or sequence specific reverse primer

Note: For random primers and oligo d(T)12-18 the final concentration in the reaction mix should be 2.5 μM .

For a sequence-specific reverse primer, the final concentration should be 200 nM.

- 10 ng to 5 μg total RNA or 1 ng to 500 ng mRNA
- 2 µl dNTP mix 20 mM total (5 mM each dATP, dGTP, dCTP and dTTP)
- RNase free water to 14 µl

EUROPE

LIEGE SCIENCE PARK • 4102 Seraing • BELGIUM • Tel.: +32 4 372 74 00 • Fax: +32 4 372 75 00 Toll-free: 00 800 666 00 123 • info@eurogentec.com • www.eurogentec.com

NORTH AMERICA

ANASPEC - 34801 Campus Drive • Fremont, CA 94555 • USA • Tel.: +1-510-791-9560 Toll-free: +1 800-452-5530 • Fax: +1 510-791-9572 • info.anaspec@deurogentec.com



Eurogentec products are sold for research or laboratory use only and are not to be administrated to humans or used for medical diagnostics.

NU-0020-50

Technical Data Sheet

- 3. Mix gently by pipeting.
- Heat mixture to 65 °C for 5 minutes in a thermocyler and quick chill on ice.
- 5. Centrifuge briefly to collect the contents to bottom of the tube and add:
 - 4 µl 5x cDNA buffer (blue cap)
 - 2 μ l RNase inhibitor at 20 U/ μ l (Optional but when using less than 100 ng of starting RNA, the addition of RNase inhibitor is essential)
- 6. Mix contents of the tube gently by pipeting and incubate 3 minutes at 37 °C in a thermocycler.
- 7. Add 0.5 1 µl (100 200 units) of Mu-MLV RT (depending on RNA amount)

Note: If less than 1 ng of RNA is used, reduce the amount of Mu-MLV RT in the reaction to 0.25 μ l (50 units) and add the RNase free water to 22 μ l final volume.

For 10 ng to 1 μ g RNA, 0.5 μ l (100 units) of Mu-MLV RT can be used

- 8. Mix gently by pipeting.
- Program the thermocycler using the following recommended parameters

Initial step¹ 10 min at 25 °C Reverse Transcriptase step 50 min at 37 °C Inactivation of the RT enzyme 5 min at 95 °C

The prepared cDNA can now be used as template for amplification in PCR.

Use only 10% of the first-strand reaction (= 2 μ l of the RT reaction) for 50 μ l PCR reaction. Adding larger amounts of the first-strand reaction may result in decreased amounts of PCR products or may also inhibit the PCR.

Related products

dNTP MIX (Lithium salts)				
Description	Quantity	Volume	Reference	
dNTP mix	1 x 20 µmoles	1 ml	NU-0010-10	
	5 x 20 µmoles	5 x 1 ml	NU-0010-50	
	10 x 20 µmoles	10 x 1 ml	NU-0010-100	
dNTP SET (Lithium salts)				
Description	Quantity	Volume	Reference	
dNTP set	4 x 25 umoles	4 x 50 ul	NU-0020-10	

5 x (4 x 25 µmoles)

Quality control

→ RNase Assay

No detectable RNase activity was observed when 10 units of the enzyme was incubated with 8 mg RNA in a 20 μ l reaction volume for 24 hours at 37 °C.

4 x 250 µl

→ Exonuclease Assay

Incubation of 50 units of the enzyme with 1 μ g Lambda DNA for 16 hours at 37 °C in the stated reaction buffer does not produce any detectable degradation of the DNA.

For further information please contact our Customer Help Desk:

For Europe:

E-mail: info@eurogentec.com Tel: +32 4 372 76 65

For USA:

E-mail: info.usa@eurogentec.com

Tel: +1 858 793 26 61

¹ Only if random primers or oligo d(T)12-18 are used.